

Item Number	Modifications for custom oligonucleotides Phosphates
MD-PH001-03004	3' Phosphate 40 nmol
MD-PH001-03020	3' Phosphate 200 nmol
MD-PH001-03100	3' Phosphate 1000 nmol
MD-PH001-05004	5' Phosphate 40 nmol
MD-PH001-05020	5' Phosphate 200 nmol
MD-PH001-05100	5' Phosphate 1000 nmol
MD-PH001-05M02	5' Phosphate 2,5 μ mol
MD-PH001-05M05	5' Phosphate 5 μ mol
MD-PH001-05M10	5' Phosphate 10 μ mol